

Batch information

MAb Anti 6His Tb cryptate Gold

HTRF MAb Anti-6HIS Tb Gold - 5K pts

Product reference :

61HI2TLA

Product characteristics :

- Lot Number	19A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	P04 100 mM pH7 + stabilizers + BSA 0.5%
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2025.01.12
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (HIS)	20.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (HIS)	5.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.00 ng
Number of tests/vial	5 000

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

